
DGK κ (DGKK)

Product code 12-408-20N

Full-length human DGK κ [1-1271(end) amino acids and D1118N of accession number NP_001013764.1] was expressed as N-terminal DYKDDDDK tagged, biotinylated protein (146 kDa) using baculovirus expression system. The protein was purified by using DYKDDDDK tag antibody agarose.

Assay platform : ADP-Glo
Substrate : Diacylglycerol

ATP (μ M) K_m app / B_{in} : 17 / 25
Metal : Mg
Reference compound : Non-disclosable
IC₅₀ at ATP B_{in} (nM) :
IC₅₀ at 1 mM ATP (nM) :
